

PROTEIN INTERACTION		
NHR	Partner Protein	Select for Profiling
AR*	AR702	
ERR α	SRCP1	
ER α *	PGC1a	
FXR*	SRCP1-2	
GR*	PGC1a	
LXR α *	SRCP1-2	
LXR β *	SRCP1-4	
MR*	PGC1a	
PPAR α *	TRAP220	
PPAR γ *	TRAP220	
PPAR δ *	TRAP220	
PR α *	SRCP1-4	
PR β *	SRCP1-4	
RAR α *	TRAP220	
RAR β *	SRCP1-2	
RXR α *	SRCP1-2	
ROR γ	SRCP 5	
RXR γ *	S4	
THR β *	TRAP220	
THR α *	SRCP1-2	
NUCLEAR TRANSLOCATION		
NHR	Cell Background	Select for Profiling
GR	CHO	
AR	HEK	
LXR α	CHO	
MR	CHO	
PR β	U2OS	

* ALSO AVAILABLE IN A SCHEDULED PANEL